

Complex IV Human Specific Activity Microplate Assay Kit ab109910

11 References [画像数 3](#)

製品の概要

製品名	Complex IV Human Specific Activity Microplate Assay Kit
検出方法	Colorimetric
サンプルの種類	Cell culture extracts, Tissue
アッセイタイプ	Enzyme activity
種交差性	交差種: Human
製品の概要	<p>This rapid multiplexing microplate kit is used to determine both the activity and quantity of cytochrome c oxidase in a human sample with greater speed and simplicity. The ratio of the two is a measure of specific activity. The COX enzyme is immunocaptured within the wells of the microplate and the activity is determined by following the oxidation of reduced Cytochrome c colorimetrically by the absorbance change at 550 nm. Subsequently, in these same wells the quantity of the enzyme can then be determined by adding a COX specific detector antibody and the appropriate alkaline phosphatase conjugated antibody; this phosphatase changes a substrate from colorless to yellow at 405 nm. This reaction takes place in a time dependent manner proportional to the amount of protein captured in the wells. Included in this kit for performance of the multiplexing assay are buffers, detergent, substrates, and 96-well microplate with monoclonal antibody pre-bound to the wells of the plate, allowing for a stream-lined assay.</p>

特記事項

Range of complex IV / cytochrome c oxidase assay kits

Biochemical assay - [ab239711](#)

Immunocapture with biochemical assay (plate-based)*** - [ab109911 \(rodent\)](#) and [ab109909 \(human\)](#)

*** Most popular assay format

Immunocapture with biochemical assay (dipstick) - [ab109878 \(rodent\)](#) and [ab109876 \(human\)](#)

Immunocapture with biochemical assay and ELISA - [ab109910 \(human\)](#) (this kit)

ELISA - [ab179880 \(human\)](#)

Other related products

Review the [mitochondrial assay guide](#), or the full [metabolism assay guide](#) to learn about more assays for metabolites, metabolic enzymes, mitochondrial function, and oxidative stress, and also how to assay metabolic function in live cells using your plate reader.

試験プラットフォーム

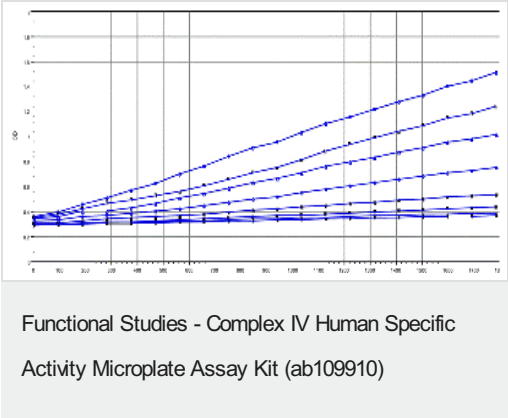
Microplate reader

製品の特性

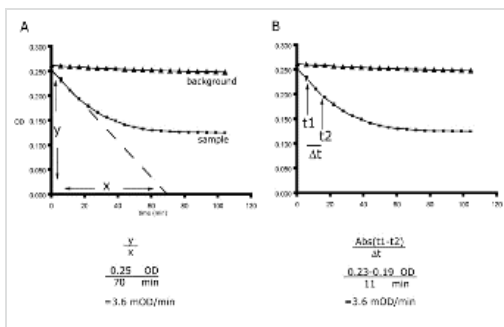
保存方法 Please refer to protocols.

内容	ラベル	96 tests
20X Buffer		1 x 15ml
96-well microplate (12 strips)		1 unit
AP Development Reagent		1 x 400µl
AP label 2500X	Tube B	1 x 12µl
Detector antibody	Tube A	1 x 1ml
Detergent		1 x 1ml
Development buffer	Tube 3	1 x 10ml
Reagent C (Reduced Cytochrome c)		1 x 1ml
Wash Buffer	Tube 2	1 x 2ml

画像

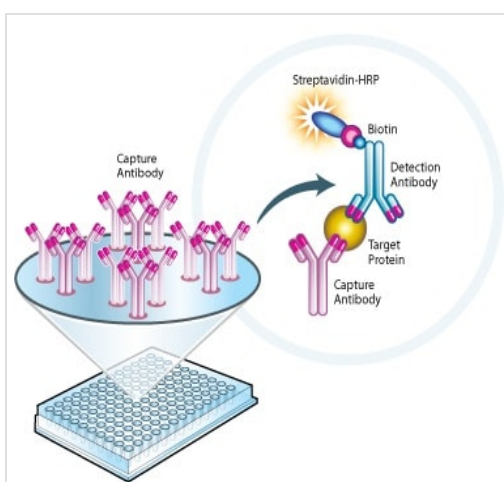


Functional study using Complex IV Human Specific Activity Microplate Assay Kit (ab109910).



Functional study using Complex IV Human Specific Activity Microplate Assay Kit (ab109910).

Functional Studies - Complex IV Human Specific Activity Microplate Assay Kit (ab109910)



Principle of Sandwich ELISA.

Sandwich ELISA - Complex IV Human Specific Activity Microplate Assay Kit (ab109910)

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.co.jp/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors